

# 20-Hydroxyecdysone

## Ecdysteroid

Most widely occurring ecdysteroid in both plant and animal species.  
Controls cell death during metamorphosis of *Drosophila melanogaster*.

Citations: 11

[View Online »](#)

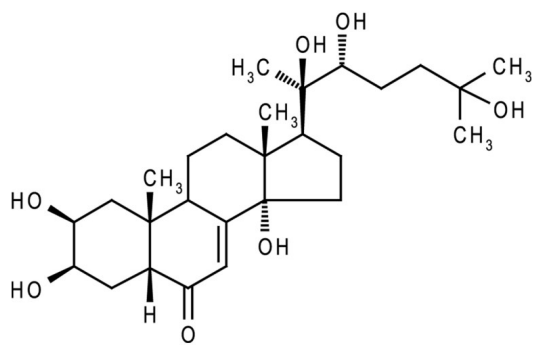
## Ordering Information

[Order Online »](#)

ALX-370-012-M005	5mg
ALX-370-012-M010	10mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name**  $\beta$ -Ecdysone, 2 $\beta$ ,3 $\beta$ ,14 $\alpha$ ,20R,22R,25-Hexahydroxy-5 $\beta$ -cholest-7-en-6-one, Ecdysterone

**Appearance** White powder.

**CAS** 5289-74-7

**Formula**  $C_{27}H_{44}O_7$

**MI** 14: 3491

**MW** 480.6

**Purity**  $\geq$ 95% HPLC

**Solubility** Soluble in water, 100% ethanol, methanol, DMSO and acetic acid.

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)